

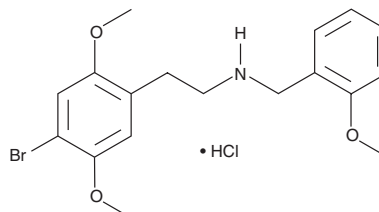
PRODUCT INFORMATION



25B-NBOMe (hydrochloride)

Item No. 12047

CAS Registry No.: 1539266-15-3
Formal Name: 4-bromo-2,5-dimethoxy-N-
[(2-methoxyphenyl)methyl]-
benzeneethanamine, monohydrochloride
Synonym: 2C-B-NBOMe
MF: C₁₈H₂₂BrNO₃ • HCl
FW: 416.7
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

2C-B is a hallucinogenic designer drug which potently activates the serotonin receptor 5-HT_{2C} (pEC₅₀ = 6.8 for arachidonic acid release).¹ 25B-NBOMe is a derivative of 2C-B, characterized by the addition of a benzyl-methoxy (BOMe) group to the amine. This N-(2-methoxybenzyl) addition significantly increases the affinity (K_i = 1.01 nM) and activity (ED₅₀ = 0.51 nM) at the serotonin receptor 5-HT_{2A}.² This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. Moya, P.R., Berg, K.A., Gutiérrez-Hernandez, M.A., *et al.* Functional selectivity of hallucinogenic phenethylamine and phenylisopropylamine derivatives at human 5-hydroxytryptamine (5-HT)_{2A} and 5-HT_{2C} receptors. *J. Pharmacol. Exp. Ther.* **321**, 1054-1061 (2007).
2. Ettrup, A., Hansen, M., Santini, M.A., *et al.* Radiosynthesis and in vivo evaluation of a series of substituted ¹¹C-phenethylamines as 5-HT_{2A} agonist PET tracers. *Eur. J. Nucl. Med. Mol. Imaging* **38**, 681-693 (2011).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/09/2019

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM